

CDNF (NT) antibody (pAb)

Rabbit Anti-Human/Mouse/Rat CDNF (NT)

Instruction Manual

Catalog Number	PK-AB718-4345
Synonyms	CDNF Antibody; Conserved dopamine neurotrophic factor, arginine-rich, mutated in early stage tumors-like 1, ARMET-like 1, ARMETL1
Description	The conserved dopamine neurotrophic factor (CDNF) is a neurotrophic factor for dopaminergic neurons and highly homologous to the mesencephalic-astrocyte-derived neuro-trophic factor. Somewhat surprisingly, CDCF is expressed at higher levels in tissues such as heart, skeletal muscle and testes than in brain. Similar to the glial cell line-derived neurotrophic factor (GDNF), CDNF can prevent the 6-hydroxylamine (6-OHDA)-induced degeneration of dopaminergic neurons in a rat model of Parkinson's disease. Furthermore, CDNF was able to restore the dopaminergic function and prevent the degeneration of dopaminergic neurons in the substantia nigra, suggesting that CDNF might be suitable for the treatment of Parkinson's disease. At least two isoforms of CDNF are known to exist. This antibody does not cross-react with MANF.
Quantity	100 µg
Source / Host	Rabbit
Immunogen	CDNF antibody was raised in rabbits against a 12 amino acid peptide from near the amino terminus of human CDNF.
Purification Method	Affinity chromatography purified via peptide column.
Clone / IgG Subtype	Polyclonal antibody
Species Reactivity	Human, Mouse, Rat
Specificity	This antibody does not cross-react with MANF.
Formulation	Antibody is supplied in PBS containing 0.02% sodium azide.
Reconstitution	During shipment, small volumes of antibody will occasionally become entrapped in the seal of the product vial. For products with volumes of 200 µl or less, we recommend gently tapping the vial on a hard surface or briefly centrifuging the vial in a tabletop centrifuge to dislodge any liquid in the container's cap.
Storage & Stability	Antibody can be stored at 4°C for three months and at -20°C for up to one year. As with all antibodies care should be taken to avoid repeated freeze thaw cycles. Antibodies should not be exposed to prolonged high temperatures.
Applications	E, WB, IHC, IF INote: Antibody might be suitable for other applications not tested so far. Optimal concentrations for each application have to be determined individually. CDNF antibody can be used for detection of CDNF by Western blot at 1 µg/mL. Antibody can also be used for immunohistochemistry starting at 5 µg/mL. For immunofluorescence start at 20 µg/mL.
Images	Available upon request.
References	Lindholm P, Voutilainen MH, Lauren J, et al. Novel neurotrophic factor CDNF protects and rescues midbrain dopamine neurons in vivo. <i>Nature</i> 2007; 448:73-7. Petrova P, Raibekas A, Pevsner J, et al. MANF: a new mesencephalic, astrocyte-derived neurotrophic factor with selectivity for dopaminergic neurons. <i>J. Mol. Neurosci.</i> 2003; 20:173-88.
Images	Available upon request.
Related Products	Cat.No. PK-AB718-4345P; CDNF Peptide Cat.No. PK-AB718-1288; A20 Cell Lysate

FOR IN VITRO RESEARCH USE ONLY. NOT FOR DIAGNOSTIC OR THERAPEUTIC PROCEDURES.